

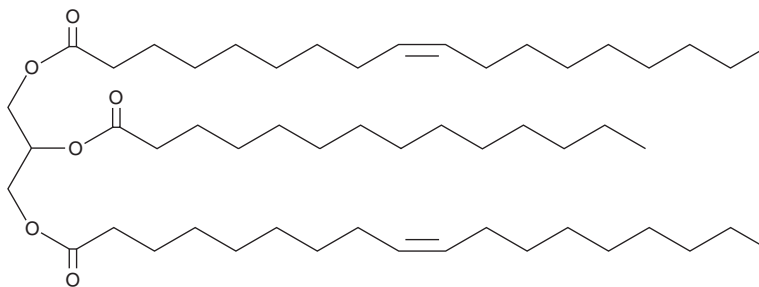
PRODUCT INFORMATION



1,3-Dioleoyl-2-Myristoyl Glycerol

Item No. 26872

CAS Registry No.: 158298-95-4
Formal Name: 9Z-octadecenoic acid, 1,1'-[2-[(1-oxotetradecyl)oxy]-1,3-propanediyl] ester
Synonyms: 1,3-Olein-2-Myristin, TG(18:1/14:0/18:1)
MF: C₅₃H₉₈O₆
FW: 831.3
Purity: ≥98%
Supplied as: An oil
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

1,3-Dioleoyl-2-myristoyl glycerol is supplied as an oil. A stock solution may be made by dissolving the 1,3-dioleoyl-2-myristoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-dioleoyl-2-myristoyl glycerol is slightly soluble in methanol.

1,3-Dioleoyl-2-myristoyl glycerol is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

1,3-Dioleoyl-2-myristoyl glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-3 positions and myristic acid (Item No. 13351) at the *sn*-2 position. It has been found in palm oil.¹

Reference

1. Kallio, H., Yli-Jokipii, K., Kurvinen, J.-P., *et al.* Regioisomerism of triacylglycerols in lard, tallow, yolk, chicken skin, palm oil, palm olein, palm stearin, and a transesterified blend of palm stearin and coconut oil analyzed by tandem mass spectrometry. *J. Agric. Food Chem.* **49**(7), 3363-3369 (2001).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/28/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM